

Fluorescein Recombinant monoclonal antibody (4-4-20 enhanced) (Human Fab fragment

This antibody is covered by our [Warranty Guarantee](#)

Manuals, SDS & CoFA

[View Online »](#)

Handling & Storage

| | |
|---------------------------|---|
| Use/Stability | Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C. |
| Handling | Avoid freeze/thaw cycles. |
| Short Term Storage | +4°C |
| Long Term Storage | -20°C |
| Shipping | Blue Ice |

Regulatory Status RUO - Research Use Only

Product Details

| | |
|---------------------------------------|--|
| Alternative Name | Fluorescyl |
| Application | Flow Cytometry, WB |
| Application Notes | Negative control |
| Clone | 4-4-20 (enhanced) |
| Formulation | Liquid. In PBS containing 0.02% Proclin 300. |
| Host | Human |
| Immunogen | Fluorescein I-bovine serum albumin (FII-BSA). |
| Isotype | Fab fragment κ |
| Purity Detail | Immobilized metal affinity chromatography purified. |
| Technical Info / Product Notes | This is a chimeric antibody created as part of a panel offering antibodies of the same specificity in different formats (species, isotype, subtype and modified versions) for use as isotype controls. |
| Worry-free Guarantee | This antibody is covered by our Worry-Free Guarantee |

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)